

Cortisol ELISA Kit ab108665

★★★★★ [4 Abreviews](#) [13 References](#) [2 Images](#)

Overview

Product name Cortisol ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall	60			<=9%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	30			<=9.8%

Sample type Urine, Serum, Hep Plasma, Cit plasma

Assay type Competitive

Sensitivity = 2.44 ng/ml

Range 10 ng/ml - 500 ng/ml

Recovery 103.56 %

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's Cortisol *in vitro* competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Cortisol in serum and plasma.

A 96-well plate has been precoated with anti-Cortisol IgG. Samples and the Cortisol-HRP conjugate are added to the wells, where any Cortisol in the sample competes with the added Cortisol-HRP for antibody binding. After incubation, the wells are washed to remove unbound material and TMB substrate is then added which is catalyzed by HRP to produce blue coloration. The reaction is terminated by addition of Stop Solution which stops the color development and produces a color change from blue to yellow. The intensity of signal is inversely proportional to the amount of Cortisol in the sample and the intensity is measured at 450 nm.

Platform Microplate

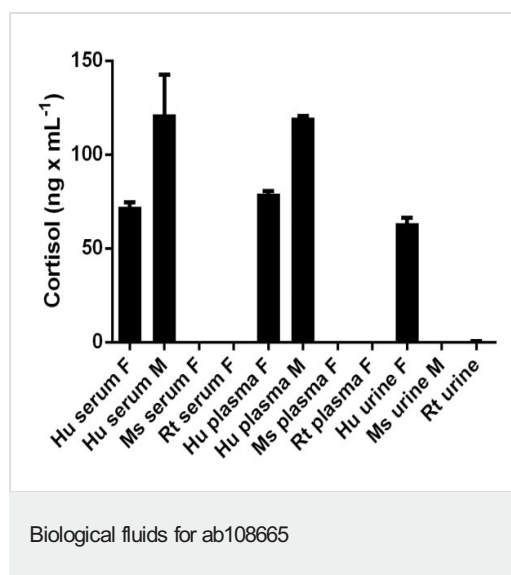
Properties

Storage instructions

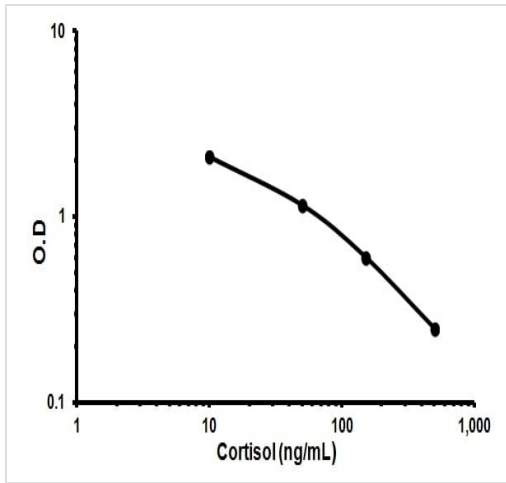
Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Washing Solution	1 x 50ml
Anti-Cortisol (IgG) Coated Microplate (12 x 8 wells)	1 unit
Cortisol Control	1 x 1ml
Cortisol Standard 0 – 0 ng/mL	1 x 1ml
Cortisol Standard 1 – 10 ng/mL	1 x 1ml
Cortisol Standard 2 – 50 ng/mL	1 x 1ml
Cortisol Standard 3 – 150 ng/mL	1 x 1ml
Cortisol Standard 4 – 500 ng/mL	1 x 1ml
Cortisol-HRP conjugate	1 x 21ml
Stop Solution	1 x 15ml
TMB Substrate Solution	1 x 15ml

Images



Cortisol measured in biological fluids showing quantity (ng) per mL of tested sample. Samples were diluted 1-2 fold.



Representative Standard Curve using ab108665

Typical Standard Curve

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors